



## Certificate of Analysis

### Mouse Interferon Beta, mammalian

**Catalog No:** 12405-1

**Lot No:** 7757

**Expiration:** July 15, 2027

**Size:**  $\geq 1 \times 10^5$  units/vial

**Description:** Recombinant Mouse Interferon Beta, mammalian

**Volume:** 0.100 ml

**Activity:**  $1.08 \times 10^6$  units/ml

**Specific Activity:**  $2.16 \times 10^8$  units/mg

**Buffer:** 20 mM Sodium Phosphate, pH 6.4; 300 mM NaCl; 5% Glycerol; 0.1% bovine serum albumin (BSA)

**Endotoxin:**  $< 1$  EU/ $\mu$ g

**Molecular Weight:** 19.6 kDa

**Purity:**  $> 95\%$

**Purification Method:** Purified from mammalian tissue culture media by a combination of ion exchange, hydrophobic interaction and size exclusion chromatography.

**Source:** Gene was obtained from mouse DNA expressed in mammalian cells

**Synonyms:** Mouse Fibroblast Interferon

**Accession #:** K00020

**Assay Used to Measure Bioactivity:** Interferon was titrated with the use of the cytopathic effect inhibition assay as listed

*Mouse (L929/EMCV)* – performed as described [Rubinstein, *et al.* (1981) *J. Virol.* 37(2):755]. The  $EC_{50}$  for interferon in this assay is  $\sim 2.5$  U/ml. Lot Activity was derived from multiple determinations in the above assay.

The units are determined by use of a Mu-IFN- $\beta$  laboratory standard calibrated to the international reference standard for mouse interferon beta provided by the National Institutes of Health [Meager, *et al.* (2001) *J. Immunol Methods.* 257 (1-2):17-33]. Please note that IFN assays vary between labs and assay systems [Meager, *et al.* (2001). *J. Immunol. Meth.* 257:17. Meager and Das (2005) *J. Immunol. Meth.* 306:1].

**Shipping Conditions:** Dry Ice

**Physical State of Product During Shipping:** Frozen

**Storage Conditions/Comments:** After receipt, the product should be kept at  $-70^{\circ}\text{C}$  or below for retention of full activity. Thaw product vial by incubation in cold tap water until just thawed – the contents of the tube should be apportioned in separate tubes so that freezing and thawing is kept to a minimum. Refreezing should be done on dry ice or in a dry ice/alcohol bath. Further dilution of the product should be in buffers containing protein such as 0.1% bovine serum albumin (BSA). For more information on protein handling, visit our Resource Library at [www.pbl assaysci.com](http://www.pbl assaysci.com).

### Authorization

Released by: \_\_\_\_\_

Date: January 31, 2025

Sold under license from Pestka Biomedical Laboratories, Inc. d/b/a PBL Assay Science. For research use only. Not for diagnostic or clinical use in, or administration to, humans. Not for resale in original or any modified form, including inclusion in a kit, for any purpose. Not for use in the preparation of any commercial product.